



IL-13 (human), ELISA kit

Catalog # **ADI-900-208**

96 Well Enzyme Immunoassay Kit

For use with serum, plasma, and cell supernatants

Table of Contents

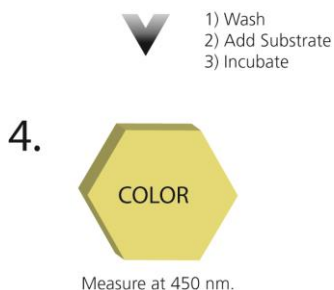
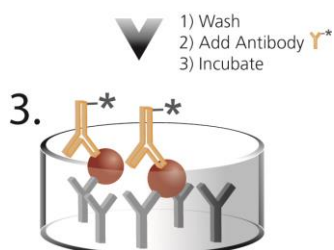
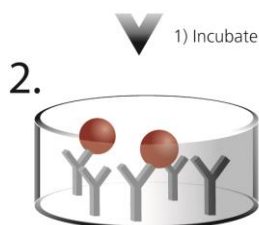
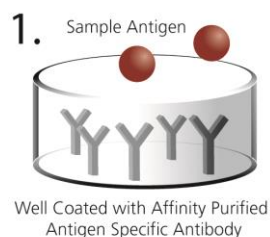


Reagents require separate storage conditions.



Check our website for additional protocols, technical notes and FAQs.

<u>2</u>	<u>Introduction</u>
<u>2</u>	<u>Principle</u>
<u>3</u>	<u>Materials Supplied</u>
<u>4</u>	<u>Storage</u>
<u>4</u>	<u>Materials Needed but Not Supplied</u>
<u>5</u>	<u>Reagent Preparation</u>
<u>6</u>	<u>Sample Handling</u>
<u>10</u>	<u>Assay Procedure</u>
<u>10</u>	<u>Calculation of Results</u>
<u>11</u>	<u>Typical Results</u>
<u>12</u>	<u>Performance Characteristics</u>
<u>13</u>	<u>References</u>
<u>16</u>	<u>Limited Warranty</u>



Introduction

The Enzo Life Sciences IL-13 Enzyme Immunometric Assay (EIA) kit is a complete kit for the quantitative determination of IL-13 in plasma, serum and cell supernatants. Please read the complete kit insert before performing this assay.

IL-13 is primarily produced by activated T cells and has many of the same biological properties as IL-4^{1,2}. IL-13 suppresses the expression of pro-inflammatory cytokines such as IL-1, IL-6, TNF- α and IL-8 by monocytes/macrophages, induces CD23 expression on B cells and promotes B cell proliferation and stimulates secretion of IgE and IgG₄ and induces IFN- γ production by NK cells^{1,2}. IL-13 has been shown to inhibit the replication of HIV both *in vitro* and *in vivo*³. IL-13 induces many phenotypes of allergic lung diseases, including airway hyper responsiveness, goblet cell metaplasia and mucus hypersecretion that contribute to airway obstruction⁴.

Principle

1. Samples and standards are added to wells coated with a monoclonal antibody to IL-13. The plate is then incubated.
2. The plate is washed, leaving only bound IL-13 on the plate. A yellow solution of horseradish peroxidase conjugated monoclonal antibody to IL-13 is then added. This binds to the IL-13 captured on the plate. The plate is then incubated.
3. The plate is washed to remove excess HRP conjugated antibody. TMB Substrate solution is added. An HRP-catalyzed reaction generates a blue color in the solution.
4. Stop solution is added to stop the substrate reaction. The resulting yellow color is read at 450nm. The amount of signal is directly proportional to the level of IL-13 in the sample.



Do not mix components from different kit lots or use reagents beyond the expiration date of the kit.



The standard should be handled with care due to the known and unknown effects of the antigen.



Protect substrate from prolonged exposure to light.



Activity of conjugate is affected by nucleophiles such as azide, cyanide, and hydroxylamine.



Stop solution is caustic. Keep tightly capped.

Materials Supplied

1. **Assay Buffer 13**
50 mL, Catalog No. 80-1500
Tris buffered saline containing BSA and detergents
2. **IL-13 Standard**
Lyophilized, Catalog No. 80-2387
2 vials containing 300 pg/vial of recombinant IL-13
3. **IL-13 Clear Microtiter Plate**
One plate of 96 wells, Catalog No. 80-2385
A clear plate of break-apart strips coated with a monoclonal antibody specific for IL-13
4. **Horseradish peroxidase IL-13 Antibody**
10ml, Catalog No. 80-2384
A yellow solution of HRP conjugated monoclonal antibody to IL-13
5. **Wash Buffer Concentrate**
27 mL, Catalog No. 80-1286
Tris buffered saline containing detergents
6. **TMB Substrate**
10 mL, Catalog No. 80-0350
A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide
7. **Stop Solution 2**
10 mL, Catalog No. 80-0377
A 1N solution of hydrochloric acid in water
8. **IL-13 Assay Layout Sheet**
1 each, Catalog No. 30-0303
9. **Plate Sealer**
3 each



Reagents require separate storage conditions.

Storage

All components of this kit, except the Standard, are stable at 4°C until the kit's expiration date. The standard must be stored at or below -20°C upon receipt.

Materials Needed but Not Supplied

1. Deionized or distilled water
2. Precision pipets for volumes between 5 μL and 1,000 μL
3. Repeater pipet for dispensing 100 μL
4. Disposable beakers
5. Graduated cylinders
6. A microplate shaker
7. Lint-free paper for blotting
8. Microplate reader capable of reading 450nm
9. Software (such as AssayBlaster™ catalog number ADI-28-0002) for extrapolating sample values from optical density readings utilizing a four parameter logistic curve fit.

Reagent Preparation



If buffers other than those provided are used in the assay, the end-user must determine the appropriate dilution and assay validation.



Standards can be made up in either glass or plastic tubes.



The standard should be handled with care due to the known and unknown effects of the antigen.

1. **Wash Buffer**

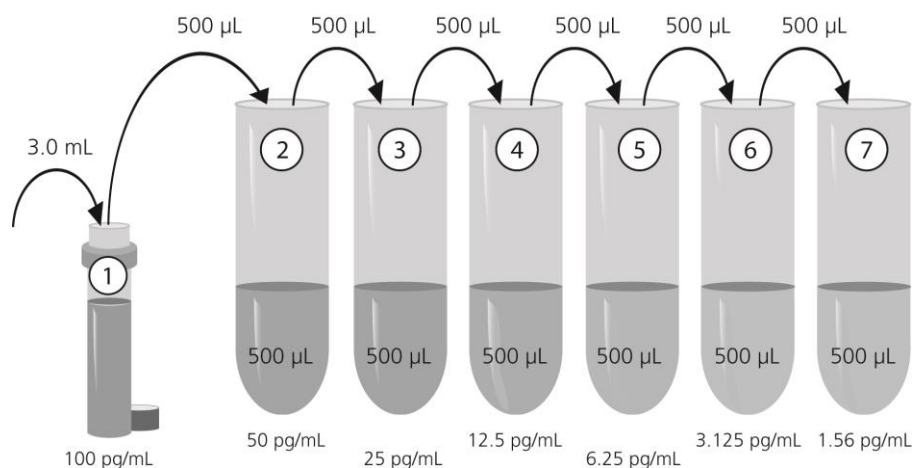
Prepare the Wash Buffer by diluting 5 mL of the supplied concentrate with 95 mL of deionized water. This can be stored at room temperature until the kit expiration date, or for 3 months, whichever is earlier.

2. **Human IL-13 Standards**

Allow the 300 pg human IL-13 standard to warm to room temperature. Reconstitute one vial of 300 pg human IL-13 Standard with 3.0ml standard diluent (Assay Buffer 13 or Tissue Culture Media) for a 100 pg/mL stock vial. Vortex thoroughly, wait 5 minutes and vortex again prior to use. Label this vial #1.

Label six disposable 12 x 75 mm tubes #2 through #7. Pipet 500 μ L standard diluent into each tube. Remove 500 μ L from reconstituted stock vial and add to tube #2 and vortex thoroughly. Remove 500 μ L from tube #2 and add to tube #3. Vortex thoroughly. Continue this for tubes #4 through #7.

The concentration of human IL-13 in tubes #1 through #7 will be 100, 50, 25, 12.5, 6.25, 3.13 and 1.56pg/ml respectively (see image below). Standards should be used within 60 minutes of preparation.



Diluted standards should be used within 1 hour of preparation. The concentrations of the standards are labeled above.

Sample Handling



Samples must be stored frozen at or below -20° to avoid loss of bio-active analyte. Repeated freeze/thaw cycles should be avoided.

The Enzo Life Sciences IL-13 (human) EIA kit is compatible with human IL-13 samples in culture supernates, plasma (with anticoagulant heparin or EDTA) and serum. Samples diluted sufficiently into the assay buffer can be read directly from a standard curve. Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens. A minimum 1:2 dilution is recommended for serum and plasma while cell culture supernates (with 10% serum supplement) may be run neat. These are the minimum recommended dilutions to remove matrix interference in the assay.

Samples in the majority of culture media, including those containing fetal bovine serum, can also be read in the assay, provided the standards have been diluted into the culture media instead of Assay Buffer. Users should only use standard curves generated in media or buffer to calculate concentrations of human IL-13 in the appropriate matrix.

Dilutional Linearity

The minimum required dilution for several common samples was determined by serially diluting samples into the assay buffer and identifying the dilution at which linearity was observed. Conditioned RPMI with 10% fetal bovine serum (FBS) was spiked with recombinant human IL-13 and diluted in Assay Buffer. The assay buffer was spiked to the same concentration and used as a control to determine linearity of the culture medium. Multiple samples of human EDTA plasma and serum were diluted in the assay buffer to produce values within the dynamic range of the assay.

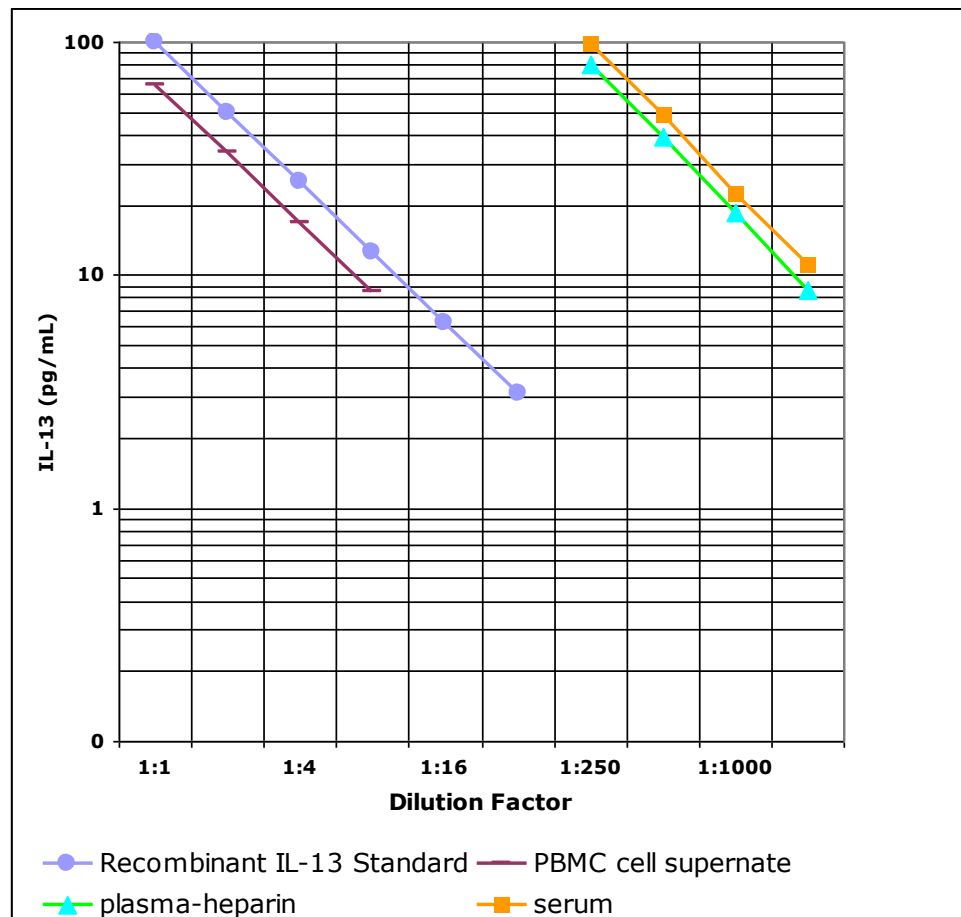
Matrix:	Media
MRD:	Neat
DF	% Dil. Linearity
Neat	93.8
1:2	96.6
1:4	98.7
1:8	92.0

Matrix:	Plasma
MRD:	1:2
DF	% Dil. Linearity
Neat	99.1
1:2	129
1:4	142
1:8	145

Matrix:	Serum
MRD:	1:2
DF	% Dil. Linearity
Neat	111
1:2	122
1:4	119
1:8	88

Parallelism

Parallelism experiments were carried out to determine if the recombinant human IL-13 standard accurately determines human IL-13 concentrations in biological matrices. Peripheral blood mononuclear cells (PBMCs) stimulated with the mitogen Phytohemagglutinin (PHA) were used to assess parallelism. Values were obtained using the cell supernatants from treated cultures serially diluted in assay buffer and assessed from a standard curve using four parameter logistic curve fitting. Human plasma and serum were similarly diluted in assay buffer and assessed again from a standard curve using four parameter logistic curve fitting. The observed values were plotted against the dilution factors. Parallelism of the curves demonstrates that the antigen binding characteristics are similar enough to allow the accurate determination of native analyte levels in diluted samples.



Spike and Recovery

After diluting each sample matrix to its minimum required dilution, recombinant human IL-13 was spiked at high, medium, and low concentrations. The recovery of the standard in spiked samples was compared to the recovery of identical spikes in the assay buffer. The mean and range of percent recovery at the three concentrations are indicated below for each matrix.

Sample Matrix	Minimum Required	Spike Concentration	Recovery of Spike
Serum (n = 2)	1:2	80	96%
		20	90%
		5	96%
Human Heparin Plasma (n=4)	1:2	80	91%
		20	86%
		5	71%
TCM + 10% FBS (n=5)	neat	80	92%
		20	85%
		10	83%

Plasma Preparation

1. Collect whole blood in vacutainer tube containing anticoagulant (heparin or EDTA).
2. Centrifuge at 1000 x g for 15 minutes at 4°C.
3. Place supernatant in a clean tube.
4. The supernatant may be divided into aliquots and stored at or below -20°C, or used immediately in the assay.
5. Avoid repeated freeze-thaw cycles.

Serum Preparation

1. Collect whole blood in appropriate tube.
2. Incubate upright at room temperature for 30-45 minutes to allow clotting to occur.
3. Centrifuge at 1000 x g for 15 minutes at room temperature. Do not use brake.
3. Without disturbing the cell layer, place supernatant into a clean tube.
4. The supernatant may be divided into aliquots and stored at or below -20°C, or used immediately in the assay.
5. Avoid repeated freeze-thaw cycles.

PBMC (Peripheral Blood Mononuclear Cells) Supernatant Preparation

1. Incubate PBMCs in RPMI supplemented with 10% FBS, L-glutamine and penicillin/streptomycin/amphotericin B for 48 hours.
2. Add phytohemagglutinin at 1 mg/mL. Incubate for 48 hours.
3. Collect cell supernatant.
4. Centrifuge at 1700 x g for 10 minutes at room temperature to pellet any cells or cellular debris.
5. Collect supernatant in a clean tube.
6. The supernatant may be divided into aliquots and stored at or below -20°C, or used immediately in the assay.
7. Avoid repeated freeze-thaw cycles.

Assay Procedure



Pre-rinse each pipet tip with reagent. Use fresh pipet tips for each sample, standard, and reagent.



Pipet the standards and samples to the bottom of the wells.



Pipet the reagents to the sides of the wells to avoid possible contamination.

Refer to the Assay Layout Sheet to determine the number of wells to be used and put any remaining wells with the desiccant back into the foil pouch and seal the ziploc. Store unused wells at 4°C.

1. Pipet 100 μL of standard diluent (Assay Buffer 13 or Culture Media) into the S0 (0pg/ml standard) wells.
2. Pipet 100 μL of Standards #1 through #7 into the appropriate wells.
3. Pipet 100 μL of the Samples into the appropriate wells.
4. Seal the plate. Incubate at room temperature on a plate shaker for 1 hour at ~500 rpm.
5. Empty the contents of the wells and wash by adding 400 μL of wash solution to every well. Repeat the wash 3 more times for a total of **4 washes**. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
6. Pipet 100 μL of yellow Antibody Solution into each well, except the Blanks.
7. Seal the plate. Incubate at room temperature on a plate shaker for 30 minutes at ~500 rpm.
8. Wash as above (Step 5). Pipet 100 μL of Substrate Solution into each well.
9. Seal the plate. Incubate at room temperature on a plate shaker for 30 minutes at ~500 rpm.
10. Pipet 100 μL of Stop Solution to each well.
11. Zero the plate reader against the Blank wells and read the optical density (OD) at 450nm, preferably with correction between 570 and 590nm. If the plate reader is not able to be blanked against the Blank wells, manually subtract the mean optical density of the Blank wells from all of the readings.

Calculation of Results



Make sure to multiply sample concentrations by the dilution factor used during sample preparation.

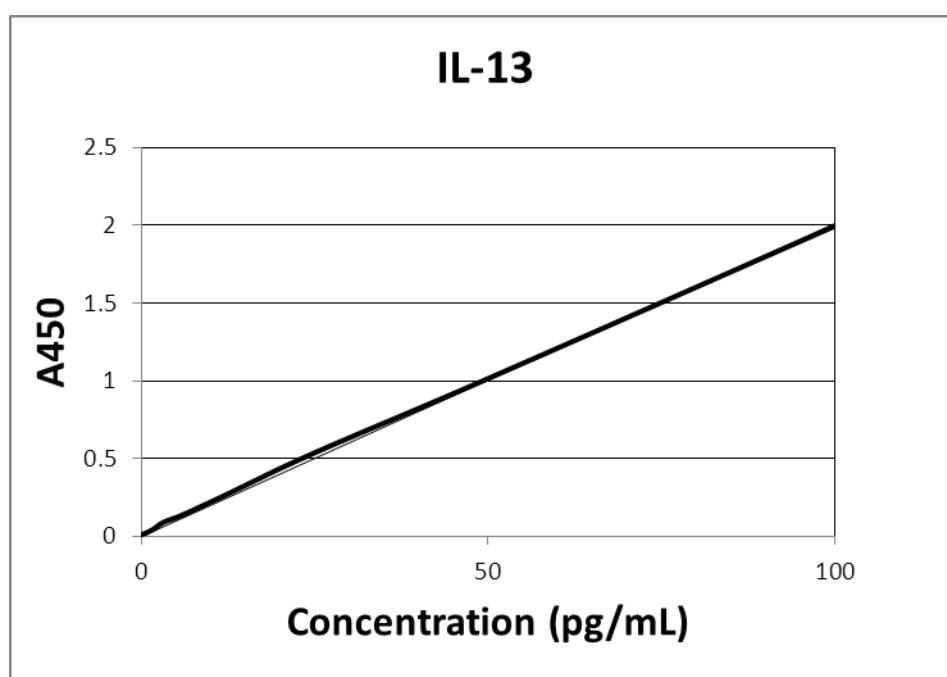
Several options are available for the calculation of the concentration of IL-13 in the samples. We recommend that the data be handled by an immunoassay software package utilizing a four parameter logistic curve fitting program (such as AssayBlaster™, catalog number ADI-28-0002). Such software is often supplied by plate reader manufacturers.

Samples with concentrations outside of the standard curve range will need to be reanalyzed using a different dilution.

Typical Results

The results shown below are for illustration only and should not be used to calculate results from another assay.

Sample	Net OD	IL-13(pg/mL)
S0	0.009	0
S1	2.0	100
S2	1.02	50
S3	0.54	25
S4	0.27	12.5
S5	0.14	6.25
S6	0.09	3.13
S7	0.04	1.56



Performance Characteristics

Specificity

The cross reactivities for a number of related compounds were determined by diluting cross reactants in the assay buffer at a concentration of 3,333-30,000 pg/mL. These samples were then measured in the assay.

Analyte	Cross Reactivity
IL-5	<0.01%
IL-15	<0.01%
IL-17 α	<0.01%
IL-21	<0.01%
IL-29	<0.01%
IL-4	<0.01%
IL-8	<0.01%
IL-1 β	<0.01%
IL-6	<0.01%
TNF α	<0.01%
RANTES	<0.01%
IFN γ	<0.01%
IL-1 α	<1.00%
IL-2	<1.00%
IL-10	<1.00%
MCP-1	<1.00%
IL-13 mouse	<1.00%

Sensitivity

The sensitivity or limit of detection of the assay is 1.71pg/ml. The sensitivity was determined by interpolation at 2 standard deviations above the mean signal at background (0pg/mL) using data from at least 6 standard curves.

Precision

Intra-assay precision was determined by assaying 20 replicates of three buffer controls containing IL-13 in a single assay.

pg/mL	%CV
51.5	6.15
12.4	2.89
2.8	7.46

Inter-assay precision was determined by measuring buffer controls of varying IL-13 concentrations in multiple assays over several days.

pg/mL	%CV
78.8	4.3
20.3	5.0
5.1	15.8

References

1. Kips, JC. Cytokines in Asthma. *European Respiratory Journal* 2001; 18: Suppl. 34, 24s-33s.
2. Elias, JA, et al. Interleukin-13 and Leukotrienes: An Intersection of Pathogenic Schema. *American Journal of Respiratory Cell and Molecular Biology* 2003; 28: 401-404.
3. Herbein, G and Varin A. The Macrophage in HIV-1 Infection: From Activation to Deactivation? *Retrovirology* 2010; 7:33 1-15.
4. Müller, U, et al. IL-13 Induces Disease-Promoting Type 2 Cytokines, Alternatively Activated Macrophages and Allergic Inflammation During Pulmonary Infection of Mice With *Cryptococcus neoformans*. *The Journal of Immunology* 2007; 179: 5367-5377.

Notes

Notes



USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

www.enzolifesciences.com
Enabling Discovery in Life Science[®]

Global Headquarters

Enzo Life Sciences Inc.
10 Executive Blvd
Farmingdale, NY 11735
(p) 1-800-942-0430
(f) 1-631-694-7501
(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lause / Switzerland
(p) +41/0 61 926 89 89
(f) +41/0 61 926 89 79

(e) info-ch@enzolifesciences.com

Please visit our website at www.enzolifesciences.com for additional contact information.